

PRODUCT INFORMATION



Voriconazole N-oxide

Item No. 28047

CAS Registry No.: 618109-05-0
Formal Name: (α R, β S)- α -(2,4-difluorophenyl)-5-fluoro- β -methyl- α -(1H-1,2,4-triazol-1-ylmethyl)-4-pyrimidineethanol, 1-oxide

Synonym: UK 121265

MF: C₁₆H₁₄F₃N₅O₂

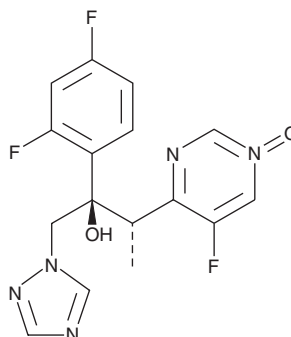
FW: 365.3

Purity: \geq 95%

Supplied as: A solid

Storage: -20°C

Stability: \geq 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Voriconazole N-oxide is supplied as a solid. A stock solution may be made by dissolving the voriconazole N-oxide in the solvent of choice, which should be purged with an inert gas. Voriconazole N-oxide is slightly soluble in chloroform and methanol.

Description

Voriconazole N-oxide is a major metabolite of the triazole antifungal voriconazole (Item No. 15633).¹ It is formed via oxidation of voriconazole by the cytochrome P450 (CYP) isoforms CYP3A4 and CYP2C19.

Reference

1. Hyland, R., Jones, B.C., and Smith, D.A. Identification of the cytochrome P450 enzymes involved in the N-oxidation of voriconazole. *Drug Metab. Dispos.* **31**(5), 540-547 (2003).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/04/2019

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM